

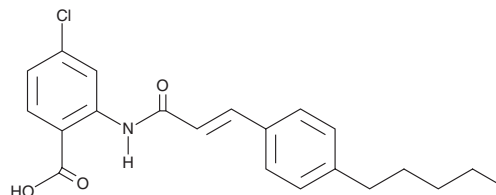
# PRODUCT INFORMATION



**ONO-RS-082**

Item No. 20243

**CAS Registry No.:** 99754-06-0  
**Formal Name:** 4-chloro-2-[[[(2E)-1-oxo-3-(4-pentylphenyl)-2-propen-1-yl]amino]-benzoic acid  
**MF:** C<sub>21</sub>H<sub>22</sub>ClNO<sub>3</sub>  
**FW:** 371.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 223, 251, 301, 321 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Special Conditions:** Light sensitive



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

## Laboratory Procedures

ONO-RS-082 is supplied as a crystalline solid. A stock solution may be made by dissolving the ONO-RS-082 in the solvent of choice, which should be purged with an inert gas. ONO-RS-082 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ONO-RS-082 in these solvents is approximately 33 mg/ml.

ONO-RS-082 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ONO-RS-082 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ONO-RS-082 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

ONO-RS-082 is a reversible inhibitor of Ca<sup>2+</sup>-independent phospholipase A<sub>2</sub>.<sup>1</sup> At 3.5 μM, it has been shown to inhibit epinephrine-stimulated thromboxane production in human platelets.<sup>1</sup> ONO-RS-082 can also disrupt endosome tubule formation and maintenance of the Golgi complex.<sup>2,3</sup>

## References

1. Banga, H. S., Simons, E. R., Brass, L. F., *et al.* Activation of phospholipases A and C in human platelets exposed to epinephrine: Role of glycoproteins IIb/IIIa and dual role of epinephrine. *Proc. Natl. Acad. Sci. USA* **83**(23), 9197-9201 (1986).
2. de Figueiredo, P., Doody, A., Polizotto, R. S., *et al.* Inhibition of transferrin recycling and endosome tubulation by phospholipase A<sub>2</sub> antagonists. *J. Biol. Chem.* **276**(50), 47361-47370 (2001).
3. Bechler, M. E. and Brown, W. J. Gβ1γ2 activates phospholipase A<sub>2</sub>-dependent Golgi membrane tubule formation. *Front. Cell. Dev. Biol.* **2**(4), 0004 (2014).

### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/14/2022

## CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM